

Agilent Array CGH

TYPE OF ARRAY	PLATFORM	SAMPLE TYPE	SAMPLE ORIGIN	REFERENCES
4x44K Agilent array CGH	Agilent Technologies	Cells	Amniotic fluid	A.L. Mosca-Boidron, et al. (2013)
105K Agilent array CGH	Agilent Technologies	Cells	Blood	A.B.P. van Kuilenburg, et al. (2010)
180K and 244K Agilent array CGH	Agilent Technologies	Cells	Blood	A. Adegbola, et al. (2015)
4x180K Agilent array CGH	Agilent Technologies	Cells	Blood	F.S. van Dijk, et al. (2010)
4x180K Agilent array CGH	Agilent Technologies	Cells	Blood	J.B. van de Meerakker, et al. (2011)
4x180K Agilent array CGH	Agilent Technologies	Cells	Blood	K. Szakson, et al. (2012)
4x180K Agilent array CGH	Agilent Technologies	Cells	Blood	E. Markljung, et al. (2012)
4x180K Agilent array CGH	Agilent Technologies	Cells	Blood	J. Aury-Landas, et al. (2013)
4x180K Agilent array CGH	Agilent Technologies	Cells	Blood	J.P. Vieira, et al. (2015)
4x44K Agilent array CGH	Agilent Technologies	Cells	Blood	B. Grisart, et al. (2008)
4x44K Agilent array CGH	Agilent Technologies	Cells	Blood	J.W. Ahn, et al. (2010)
4x44K Agilent array CGH	Agilent Technologies	Cells	Blood	A. Hills, et al. (2010)
4x44K Agilent array CGH	Agilent Technologies	Cells	Blood	S. Hoang, et al. (2011)
60K Agilent array CGH	Agilent Technologies	Cells	Blood	M. Tropeano, et al. (2013)
60K Agilent array CGH	Agilent Technologies	Cells	Blood	J.W. Ahn, et al. (2014)
4x44K Agilent array CGH	Agilent Technologies	Cells	Gastric cancer cell lines	T.E. Buffart, et al. (2009)
4x44K Agilent array CGH	Agilent Technologies	Cells	Huntington's disease human embryonic stem cells	L. Jacquet, et al. (2015)
60K Agilent array CGH	Agilent Technologies	Cells	Induced pluripotent stem cells	T. Yagi, et al. (2011)
180K and 244K Agilent array CGH	Agilent Technologies	Cells	Lymphoblastoid cell lines	A. Adegbola, et al. (2015)
60K Agilent array CGH	Agilent Technologies	ChIP DNA	ChIP-seq DNA library	D. Schmidt, et al. (2008)
4x180K Agilent array CGH	Agilent Technologies	FFPE tissue	Adenoid cystic carcinoma	A.F. Costa, et al. (2014)
4x180K Agilent array CGH	Agilent Technologies	FFPE tissue	Epithelial ovarian tumors	M.M. Kamieniak, et al. (2013)
4x44K Agilent array CGH	Agilent Technologies	FFPE tissue	Polypoid and flat colorectal adenoma	Q.J. Voorham, et al. (2012)
4x44K Agilent array CGH	Agilent Technologies	FFPE tissue	Stage II colon cancer samples	R.P. Brosens, et al. (2011)

Complete citation can be found on product's webpage

Advance Your Research:
www.enzolifesciences.com

GLOBAL HEADQUARTERS

Enzo Life Sciences Inc.
10 Executive Boulevard
Farmingdale, NY 11735
Phone: 800.942.0430
Fax: 631.694.7501
info-usa@enzolifesciences.com

EUROPE

Enzo Life Sciences (ELS) AG
Industriestrasse 17
CH-4415 Lausen, Switzerland
Phone: +41 61 926 8989
Fax: +41 61 926 8979
info-eu@enzolifesciences.com

NimbleGen & Oxford Array CGH

TYPE OF ARRAY	PLATFORM	SAMPLE TYPE	SAMPLE ORIGIN	REFERENCES
180K Oxford array CGH	Agilent Technologies	Cells	Blood	J. Winberg, et al. (2014)
1M Oxford array CGH	Agilent Technologies	Cells	Blood	A. Norling, et al. (2013)
2x105K Oxford array CGH	Agilent Technologies	Cells	Blood	D.Q. Barge-Schaapveld, et al. (2011)
135K NimbleGen array CGH	Agilent Technologies	FFPE tissue	Breast cancer tissue	P.C. Schouten, et al. (2015)
2x105K Oxford array CGH	Agilent Technologies	FFPE tissue	Primary ovarian tumor tissue	A. Zaal, et al. (2012)
180K Oxford array CGH	Agilent Technologies	Frozen tissue	Heart, liver, lung and spleen tissue	J. Winberg, et al. (2014)

BAC Array CGH

TYPE OF ARRAY	PLATFORM	SAMPLE TYPE	SAMPLE ORIGIN	REFERENCES
BAC array CGH	Agilent Technologies	Cells	Blood	S. Struski, et al. (2007)
BAC array CGH	Molecular Devices	Cells	Neuroblastoma cell lines	H. Auer, et al. (2007)
BAC array CGH	Molecular Devices	FFPE tissue	Melanoma and melanocytic spitz nevi	L. Ali, et al. (2010)
BAC array CGH	Molecular Devices	FFPE tissue	Microdissected intraductal and invasive breast cancer	S. Liao, et al. (2012)
BAC array CGH	Molecular Devices	Fresh tissue	Mouse tail	Z. Li, et al. (2007)
BAC array CGH	Molecular Devices	Frozen tissue	Primary uveal Melanoma	L.A. Worley, et al. (2007)

Other Array Types

TYPE OF ARRAY	PLATFORM	SAMPLE TYPE	SAMPLE ORIGIN	REFERENCES
Exon-level array CGH	Agilent Technologies	Cells	Blood	S. Aradhya, et al. (2012)
Large-insert clone whole-genome tilepath array	Agilent Technologies	Cells	Sperm	D.J. Tuner, et al. (2008)
<i>D.melanogaster</i> NimbleGen array CGH	NimbleGen	ChIP DNA	Drosophila S2 cells	H. Mason-Suares, et al. (2013)
<i>P.gingivalis</i> strain W83 array CGH	Agilent Technologies	Frozen cells	<i>P.gingivalis</i>	J. Brunner, et al. (2010)